

Analyzing Vapor Phase Samples with the Nicolet iS50 Multi Range Spectrometer

Key Words

ABX Automated Beamsplitter Exchanger, Gas Phase Research, Multi-spectral Range, Quality Control Applications

Fourier transform spectroscopy was initially developed by molecular physicists to provide a way to observe the high resolution structure in the infrared spectral region of gas phase molecules. Instruments were designed to measure not only chemicals in the atmosphere but also at infrared emissions throughout the universe. Infrared spectroscopy has been used extensively in physical chemistry to characterize gas phase molecules and as a routine analytical technique to determine the concentrations of specific compounds in commercial products and processes, as well as environmental monitoring. A major benefit of infrared spectroscopy is that both oxygen and nitrogen are transparent in the mid-infrared region, resulting in very low detection limits for trace chemicals that are infrared active. However, both water and carbon dioxide can be interferences, particularly in combustion analysis, where they are major reaction products. In many applications, infrared spectroscopy can routinely detect parts per billion (ppb) levels of specific chemicals.

In this application note, several examples are described where the Thermo Scientific Nicolet iS50 spectrometer has proven valuable in characterizing gas phase samples. Major areas of interest are:

- Physical chemistry and molecular characterization
- Quality assurance/quality control (QA/QC) and contaminant detection
- Combustion and catalyst research
- Environmental and air monitoring

We will also describe spectra showing the expanded spectral range capability of the Nicolet™ iS™50 spectrometer into the far-infrared and visible regions of the spectrum.



Figure 1

Experimental

All spectra reported in this paper were acquired on a Nicolet iS50 spectrometer using various pathlength gas cells from 5 centimeters to 10 meters. Spectral resolution and scan times were optimized for each application. The specific instrument used in this work was configured with all-aluminum optics for operation in the visible spectral region and the iS50 ABX automated beamsplitter exchanger containing KBr ($7000\text{--}400\text{ cm}^{-1}$), solid substrate ($700\text{--}100\text{ cm}^{-1}$) and Quartz ($25,000\text{--}5000\text{ cm}^{-1}$) beamsplitters. The combination of the three beamsplitters, the three software-selectable detectors (visible, mid-IR and far-IR) and two sources (visible and infrared) provides a fully integrated, high resolution spectrometer that can automatically cover the spectrum from $25,000\text{--}100\text{ cm}^{-1}$. Figure 1 shows the Nicolet iS50 spectrometer with 10 meter gas cell sample compartment however many different gas cells can be precisely mounted in the sample compartment.

Some users dedicate an instrument to gas analysis while most users in analytical, or QC laboratories, deal with multiple samples including liquids and solids. One of the greatest benefits of the Nicolet iS50 is flexibility and the ability to rapidly switch between sampling options. Figure 1 also shows a Nicolet iS50 configured with the ABX automatic beamsplitter exchanger which provides the rapid and precise switching of the optical components required for different spectral ranges. The 10 meter gas cell can be mounted in the sample compartment and easily swapped with other accessories including the iS50 Raman module. All Nicolet iS50 spectrometer systems are validated with a CO sample during final test. Figure 2 shows a spectrum acquired with the 5 cm CO standard cell with a full width at half height of better than 0.1 cm^{-1} .

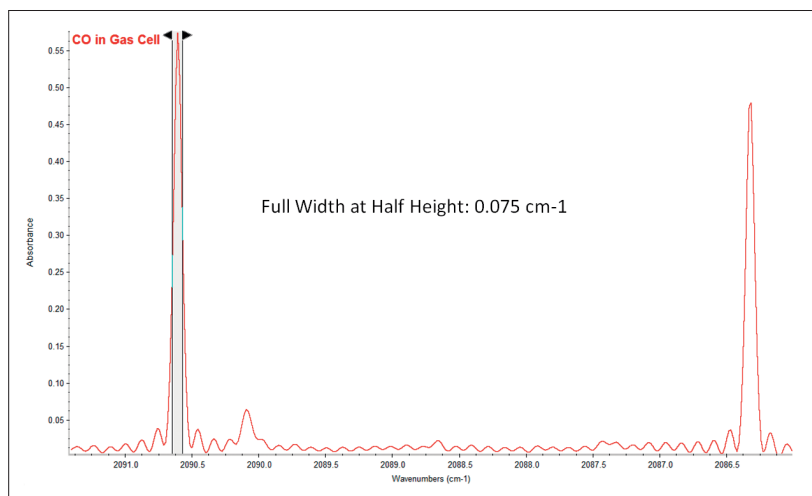


Figure 2

Physical Chemistry and Molecular Studies

The infrared spectra of gas phase molecules are often a very complex set of peaks corresponding not only to vibrational transitions but also the quantized rotational transitions of the molecule. These peaks often provide valuable information about the chemical bonds and structure of the molecule. This unique combination of vibrational and rotational transitions also provides the spectral specificity that makes multicomponent gas analysis possible. A good example that demonstrates the capability of the Nicolet iS50 spectrometer is a traditional physical chemistry laboratory experiment involving the visible spectrum of iodine vapor. While iodine is a solid at room temperature, it easily sublimates and an excellent spectrum can be obtained by placing iodine crystals in the bottom of a 10 cm gas cell. Although iodine has no allowed infrared vibrational modes, it has a multitude of spin-forbidden transitions between the vibrational levels of the ground state and levels in the excited triplet states. The result is a unique high resolution spectrum from $20,000\text{--}17,000\text{ cm}^{-1}$ as shown in Figure 3.

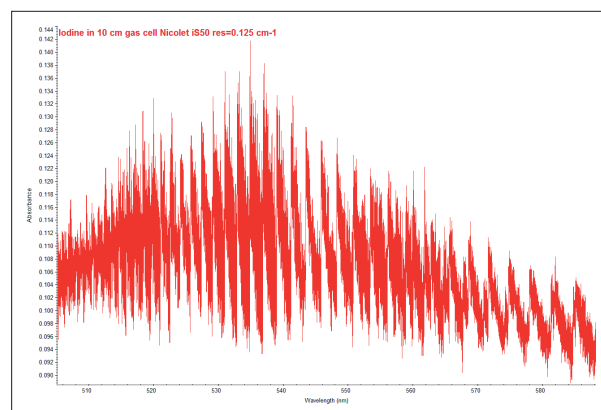


Figure 3

A second example used in physical chemistry is the mid-infrared spectrum of ammonia. The spectrum was acquired in a 10 cm cell at 0.125 cm^{-1} spectral resolution with the Nicolet iS50 as shown in Figure 4. The doubling of the Q-branch due to inversion is clearly seen, as are the multiple high resolution bands for rotations.

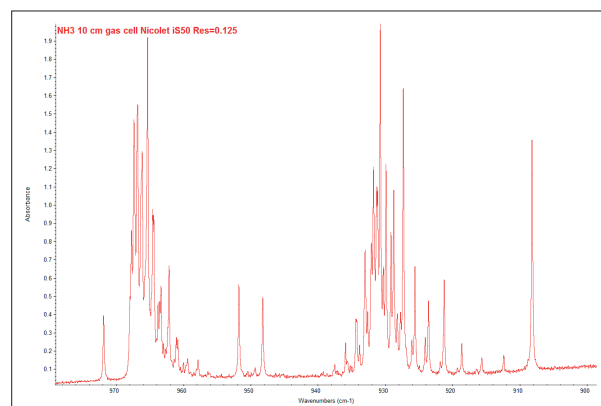


Figure 4

Far-infrared spectroscopy has also been used extensively to characterize gas phase molecules. The high-resolution spectrum shown in Figure 5 was collected in a single-button-push on the Nicolet iS50, switching the beamsplitter from the visible spectral region to the far-infrared.

QA/QC and Air Purity

A very important U.S. QA/QC application of infrared gas analysis is screening for trace contamination in Aviator's Breathing Oxygen (ABO). Both the Air Force and the Navy have systems stationed throughout the world to verify that no dangerous contaminants are present in the oxygen. Figure 6 lists the 20 components analyzed with this method.

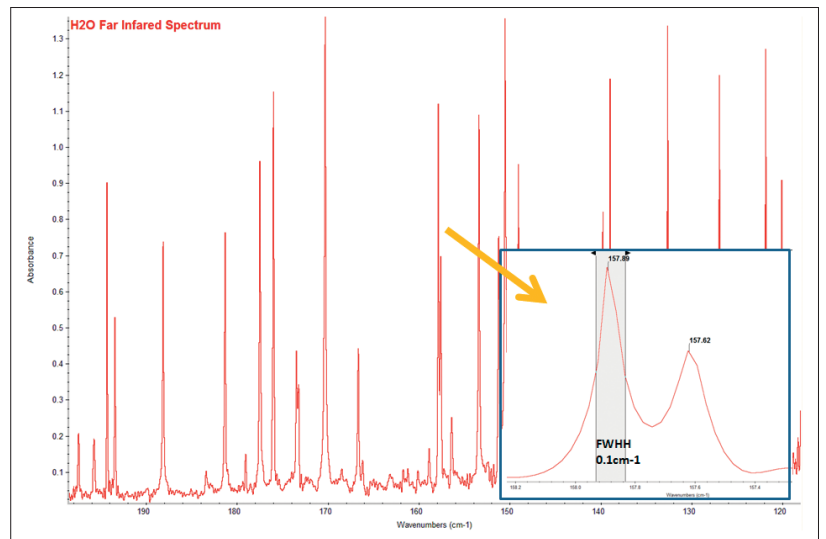


Figure 5

Aviator's Breathing Oxygen Method and Calibration

The Aviator's Breathing Oxygen (ABO) method is designed to detect impurities in ABO gas according to the U.S. Air Force military standard 1564A. This method is used with the 10 meter gas cell. Detection limits assume a collection time of 2 minutes with a room-temperature DTGS detector.



Figure 6

| | Type of Calibration | Calibration Range | Notes |
|-----------------------|---------------------------------------|-------------------|-------|
| Acetylene | One point calibration from a cylinder | 0.5 – 10 ppm | |
| Carbon dioxide | One point calibration from a cylinder | 1 – 20 ppm | |
| Carbon monoxide | One point calibration from a cylinder | 0.1 – 10 ppm | |
| CFC 11 | One point calibration from a cylinder | 0.5 – 10 ppm | |
| CFC 113 | One point calibration from a cylinder | 0.5 – 10 ppm | |
| CFC 12 | One point calibration from a cylinder | 0.5 – 10 ppm | |
| CFC 13 | One point calibration from a cylinder | 0.5 – 10 ppm | |
| CFC 141b | One point calibration from a cylinder | 0.5 – 10 ppm | |
| CFC 22 | One point calibration from a cylinder | 0.5 – 10 ppm | |
| CFC 225 | One point calibration from a cylinder | 0.5 – 10 ppm | |
| Ethane | One point calibration from a cylinder | 0.5 – 10 ppm | |
| Ethylene | One point calibration from a cylinder | 0.5 – 20 ppm | |
| Methane | One point calibration from a cylinder | 0.5 – 50 ppm | |
| Nitrous oxide | One point calibration from a cylinder | 0.5 – 10 ppm | |
| Propane | One point calibration from a cylinder | 0.5 – 10 ppm | |
| Propylene | One point calibration from a cylinder | 0.5 – 10 ppm | |
| Sulfur hexafluoride | One point calibration from a cylinder | 0.5 – 10 ppm | |
| 1,1,1-Trichloroethane | One point calibration from a cylinder | 0.5 – 10 ppm | |
| Trichloroethylene | One point calibration from a cylinder | 0.5 – 10 ppm | |
| Water | Factory calibrated (Room temp – 60°C) | 5 – 100 ppm | 1 |

NOTES: 1. Use ZnSe if greater than 3% water. Do not use ZnSe if more than a few hundred ppm of halogens (e.g.: HF, HCl, HBr) or % level acidic gases, such as NO.

In this example, spectra were acquired in a 10 meter gas cell on the Nicolet iS50 from gas standards and verification samples used to calibrate the Air Force method. Spectra were acquired from each cylinder with a one minute scan time at 1 cm⁻¹ resolution as shown in Figure 7. A validation sample containing 10 ppm levels of these components was provided by the U.S. Navy as part of this project.

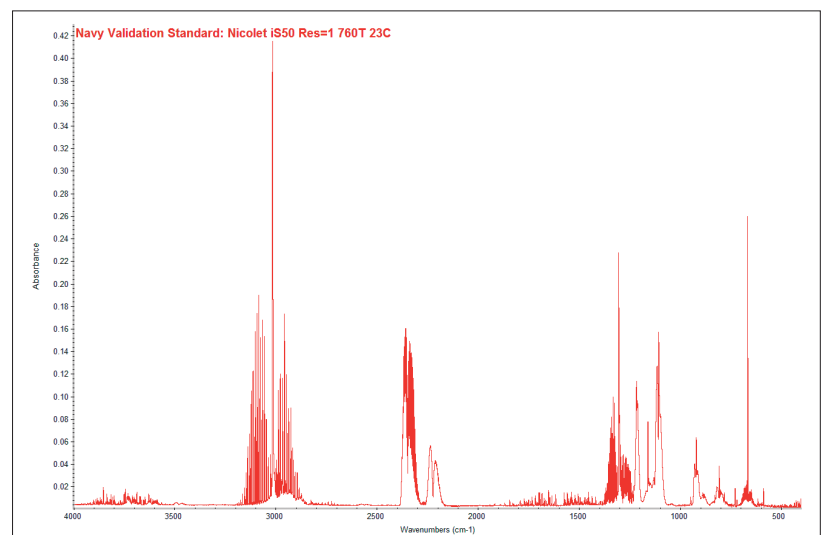


Figure 7

A Thermo Scientific TQ Analyst software method was calibrated with spectra obtained from gas standards purchased from Scott Specialties Gases® and tested on a spectrum of the Navy test gas. The method was also applied to a spectrum obtained from a nitrogen zero gas. The results are shown in Figure 8.

| Navy Reference Gas Mixture | | | | Nitrogen Zero Gas | | | |
|----------------------------|-----------------------|---------------|------------|-------------------|-----------------------|---------------|------------|
| Index | Component | Concentration | Std. error | Index | Component | Concentration | Std. error |
| 1 | Methane | 48.62 | 0.224 | 1 | Methane | -0.01 | 0.006 |
| 2 | Ethane | 5.97 | 0.320 | 2 | Ethane | -0.02 | 0.011 |
| 3 | Ethylene | 0.82 | 0.043 | 3 | Ethylene | 0.03 | 0.033 |
| 4 | Acetylene | 0.24 | 0.022 | 4 | Acetylene | 0.01 | 0.015 |
| 5 | Propane | 0.00 | 0.208 | 5 | Propane | 0.00 | 0.007 |
| 6 | Propene | 0.00 | 0.094 | 6 | Propene | -0.05 | 0.065 |
| 7 | Freon-11 | 0.03 | 0.007 | 7 | Freon-11 | 0.01 | 0.005 |
| 8 | Freon-12 | 2.26 | 0.058 | 8 | Freon-12 | -0.01 | 0.006 |
| 9 | Freon-13 | 1.98 | 0.012 | 9 | Freon-13 | -0.00 | 0.001 |
| 10 | Freon-22 | 2.84 | 0.025 | 10 | Freon-22 | 0.03 | 0.020 |
| 11 | Freon-113 | 0.39 | 0.073 | 11 | Freon-113 | -0.00 | 0.008 |
| 12 | 1,1,1-Trichloroethane | 0.00 | 0.140 | 12 | 1,1,1-Trichloroethane | 0.01 | 0.097 |
| 13 | Tetrachloroethylene | 0.00 | 0.031 | 13 | Tetrachloroethylene | -0.00 | 0.024 |
| 14 | Nitrous oxide | 4.16 | 0.033 | 14 | Nitrous oxide | 0.00 | 0.005 |
| 15 | Carbon monoxide | 0.00 | 0.066 | 15 | Carbon monoxide | 0.20 | 0.049 |
| 16 | Carbon dioxide Low | 10.28 | 0.016 | 16 | Carbon dioxide Low | 0.06 | 0.008 |
| 17 | Water | 6.65 | 0.073 | 17 | Water | 0.48 | 0.057 |
| 18 | Freon-141b | 0.00 | 0.122 | 18 | Freon-141b | 0.02 | 0.014 |
| 19 | Sulfur hexafluoride | 0.00 | 0.001 | 19 | Sulfur hexafluoride | -0.00 | 0.001 |
| 20 | Freon-225 | 0.00 | 0.111 | 20 | Freon-225 | 0.00 | 0.013 |

Figure 8

These results are reported as ppm and show a standard error of less than 100 ppb for most of the 20 components. The standard errors are even lower for the results with zero gas with the calculated values for all components below 1 ppm and many below 100 ppb.

Similar methods have been used in many industries to detect trace contamination and moisture. Many producers of specialty gases use infrared spectroscopy as a key analysis technique in their QA/QC laboratories.

Environmental and Air Monitoring

Infrared spectroscopy has been used extensively to detect volatile organic species and other pollutants in the air. These hazardous compounds may result from a manufacturing process, land fill off-gas, vehicle exhaust or a chemical spill. A major challenge in measuring pollutants in the air is the strong infrared absorbance for both water and CO₂ which block out many of the spectral regions. A spectral resolution of 0.5 cm⁻¹ is often necessary to create analysis windows between the water and CO₂ peaks as shown in Figure 9.

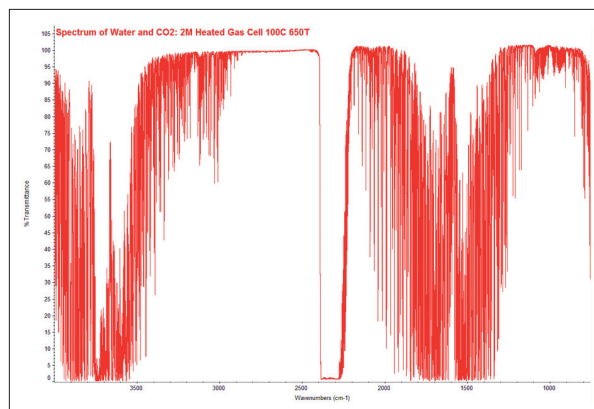


Figure 9

In many applications, continuous extractive sampling, Summa canisters or Tedlar bags are used to collect samples which can be pulled into the same 10 meter gas cell described previously. In high humidity situations the gas cell may be heated above 100 °C to ensure that water does not condense in the gas cell.

Conclusion

In this article, we have presented an overview of various gas analysis application areas where infrared spectroscopy has proven valuable. The Nicolet iS50 is designed with flexibility to analyze a variety of sample types with ease. The combination of high sensitivity and a full suite of software features, specifically designed for gas analysis, creates a world class solution to a broad range of applications.

thermoscientific.com

©2012 Thermo Fisher Scientific Inc. All rights reserved. Scott Gas is a part of Air Liquide America Specialty Gases LLC. Tedlar is a registered trademark of E. I. du Pont de Nemours and Company. ISO is a trademark of the International Standards Organization. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.



Thermo Electron Scientific Instruments LLC, Madison, WI
USA is ISO Certified.

| | | | |
|----------------------------------|--|--------------------------------------|-------------------------------------|
| Africa +27 11 822 4120 | Denmark +45 70 23 62 60 | India +91 22 6742 9434 | New Zealand +64 9 980 6700 |
| Australia +61 3 9757 4300 | Europe-Other +43 1 333 50 34 0 | Italy +39 02 950 591 | Russia/CIS +43 1 333 50 34 0 |
| Austria +43 1 333 50 34 0 | Finland/Norway/Sweden +46 8 556 468 00 | Japan +81 45 453 9100 | Spain +34 914 845 965 |
| Belgium +32 53 73 42 41 | France +33 1 60 92 48 00 | Latin America +1 561 688 8700 | Switzerland +41 61 716 77 00 |
| Canada +1 800 530 8447 | Germany +49 6103 408 1014 | Middle East +43 1 333 50 34 0 | UK +44 1442 233555 |
| China +86 10 8419 3588 | | Netherlands +31 76 579 55 55 | USA +1 800 532 4752 |

AN52316_E 04/12M

Thermo
SCIENTIFIC

Part of Thermo Fisher Scientific